

PRODUCT INFORMATION



Proadrenomedullin (N-20) (rat) (trifluoroacetate salt)

Item No. 27405

Formal Name:	L-alanyl-L-arginyl-L-leucyl-L- α -aspartyl-L-threonyl-L-seryl-L-seryl-L-glutaminy-L-phenylalanyl-L-arginyl-L-lysyl-L-lysyl-L-tryptophyl-L-asparaginy-L-lysyl-L-tryptophyl-L-alanyl-L-leucyl-L-seryl-L-argininamide, trifluoroacetate salt
Synonyms:	ProADM, ProAM, Mid-Regional proADM, Mid-Regional Proadrenomedullin, MR-proADM, MR-proAM H—Ala—Arg—Leu—Asp—Thr—Ser—Ser—Gln—Phe—Arg—Lys—Lys—Trp—Asn—Lys—Trp—Ala—Leu—Ser—Arg—NH ₂ • XCF ₃ COOH
MF:	C ₁₁₁ H ₁₇₇ N ₃₇ O ₂₈ • XCF ₃ COOH
FW:	2,477.8
Purity:	≥95%
Supplied as:	A solid
Storage:	-20°C
Stability:	≥4 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Proadrenomedullin (N-20) (rat) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the proadrenomedullin (N-20) (rat) (trifluoroacetate salt) in water. The solubility of proadrenomedullin (N-20) (rat) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Proadrenomedullin (N-20) is a peptide comprised of amino acid residues 22-41 of rat proadrenomedullin.¹ The rat proadrenomedullin (N-20) sequence is similar to human proadrenomedullin (N-20) (Item No. 27102) but contains three amino acid substitutions at residues 5, 6, and 8 of the human proadrenomedullin (N-20) sequence.

Reference

1. Sakata, J., Shimokubo, T., Kitamura, K., *et al.* Molecular cloning and biological activities of rat adrenomedullin, a hypotensive peptide. *Biochem. Biophys. Res. Commun.* **195**(2), 921-927 (1993).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/20/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM